

CERTIFICATE OF ANALYSIS

Page 1 of 3

Product Description: Muscle Injury Panel 3 (mouse) Kit
Kit Catalog Numbers: K15186C-Series
Kit Lot Number: K0040442
Expiration Date: 31 Mar 2019

Acceptance Data/ Percent Recovery Specifications:

Spot	Name	Calibrator		Muscle Injury Panel 3 (mouse) Control 1		Muscle Injury Panel 3 (mouse) Control 2		Muscle Injury Panel 3 (mouse) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	cTnI	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
3	FABP3	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
5	MyI3	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%
6	sTnI	80% - 120%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%	75% - 125%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	cTnI	0.0160	75% - 125%	< 25%	20.0	80% - 120%	< 20%
3	FABP3	0.535	80% - 120%	< 20%	26.8	80% - 120%	< 20%
5	MyI3	0.135	80% - 120%	< 20%	41.4	80% - 120%	< 20%
6	sTnI	0.505	80% - 120%	< 20%	155	80% - 120%	< 20%

Acceptance Data/ Non Specific Binding Specifications:

Spot	Name	NSB Spec
1	cTnI	< 1.0%
3	FABP3	< 1.0%
5	MyI3	< 1.0%
6	sTnI	< 1.0%

CERTIFICATE OF ANALYSIS

Page 2 of 3

Stock Concentration of Material Tested :

Spot	Name	Item	Conc	Units
1	cTnI	20X Calibrator	480	ng/mL
		Muscle Injury Panel 3 (mouse) Control 1	43.2	ng/mL
		Muscle Injury Panel 3 (mouse) Control 2	5.88	ng/mL
		Muscle Injury Panel 3 (mouse) Control 3	1.6	ng/mL
3	FABP3	20X Calibrator	1966	ng/mL
		Muscle Injury Panel 3 (mouse) Control 1	48.6	ng/mL
		Muscle Injury Panel 3 (mouse) Control 2	12.2	ng/mL
		Muscle Injury Panel 3 (mouse) Control 3	4.64	ng/mL
5	MyI3	20X Calibrator	1300	ng/mL
		Muscle Injury Panel 3 (mouse) Control 1	105	ng/mL
		Muscle Injury Panel 3 (mouse) Control 2	26.6	ng/mL
		Muscle Injury Panel 3 (mouse) Control 3	1.8	ng/mL
6	sTnI	20X Calibrator	3860	ng/mL
		Muscle Injury Panel 3 (mouse) Control 1	195	ng/mL
		Muscle Injury Panel 3 (mouse) Control 2	31.6	ng/mL
		Muscle Injury Panel 3 (mouse) Control 3	5.04	ng/mL

Controls tested may not be the same lot as shipped.

Concentrations provided in this table refer to the stock concentration of the calibrators. For top of the standard curve concentrations, please refer to the Standard Curve Data

Standard Curve Data:

cTnI - Spot 1			FABP3 - Spot 3			MyI3 - Spot 5		
Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV	Conc. ng/mL	ECL Signal	%CV
24.0	554,000	4.4	98.3	23,600	5.0	65.0	109,000	5.7
6.00	120,000	7.0	24.6	16,400	3.9	16.3	28,800	9.0
1.50	25,400	5.2	6.14	5,690	2.4	4.06	6,640	6.6
0.375	5,720	6.8	1.54	921	1.8	1.02	1,630	11.0
0.0938	1,490	6.4	0.384	354	4.4	0.254	549	5.3
0.0234	456	5.8	0.0960	298	6.8	0.0635	258	5.4
0.00586	197	4.7	0.0240	300	7.9	0.0159	205	4.9
0.00	113	12.0	0.00	327	9.8	0.00	175	12.0

CERTIFICATE OF ANALYSIS

Page 3 of 3


sTnI - Spot 6		
Conc. ng/mL	ECL Signal	%CV
193	686,000	4.4
48.3	186,000	3.3
12.1	42,700	2.5
3.02	10,100	2.0
0.754	2,940	4.0
0.188	961	5.9
0.0471	510	9.3
0.00	289	11.0

Components:

Description	Lot Number	Storage Temperature	Expiration Date
Muscle Injury Panel 3 (Mouse) Plate	Z0055960	2-8°C	30 Nov 2019
Anti-ms cTnI Antibody	D0040712	2-8°C	30 Nov 2019
Anti-ms FABP3 Anibody	D0040643	2-8°C	31 Mar 2019
Anti-ms Myl3 Antibody	D0040714	2-8°C	31 Jan 2020
Anti-ms sTnI Antibody	D0040715	2-8°C	30 Nov 2019
Muscle Injury Panel 3 (mouse) Calibrator Blend	A0040196	≤ -70°C	30 Nov 2019
Diluent 33	M0330010	≤ -10°C	31 Dec 2019
Diluent 30	M0300021	≤ -10°C	30 Jun 2020
DTT	R0010598	≤ -10°C	31 Dec 2019
EDTA pH 8.0	R00X1122	Room Temperature	31 Mar 2020
MSD Read Buffer T	Not Kit Specific	Not Kit Specific	N/A

Statement:

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance		21 AUG 2018